

Mouse Laminin ELISA Kit ab119572

6 References 1 图像

概述

产品名称	小鼠Laminin ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Sandwich (quantitative)
灵敏度	< 10 pg/ml
范围	156 pg/ml - 10000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	<p>Abcam's mouse Laminin <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse Laminin in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).</p> <p>A polyclonal antibody from rabbit specific for Laminin has been precoated onto 96-well plates. Standards(from murine sarcoma basement membrane) and test samples are added to the wells, a biotinylated detection polyclonal antibody from rabbit specific for Laminin is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the mouse Laminin amount of sample captured in plate.</p>
平台	Microplate

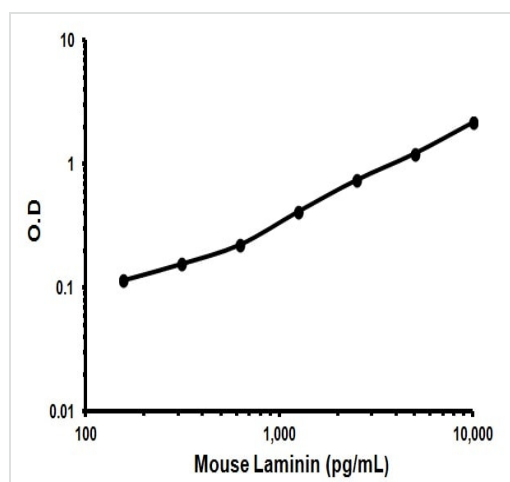
性能

存放说明Store at -20°C. Please refer to protocols.

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml

组件	标识符	1 x 96 tests	1 x 96 tests
Anti-mouse Laminin Antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100μl	1 x 100μl
Biotinylated anti-mouse Laminin antibody		1 x 100μl	1 x 100μl
Lyophilized Recombinant mouse Laminin standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

图片



Typical Standard Curve

Representative Standard Curve using ab119572.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors